



Product Information Sheet

Product Name:	Goat Anti-Mouse IgG (H+L), highly cross-adsorbed, Hilyte Fluor™ 488-labeled
Catalog Number:	61057-05-H488
Size:	0.5 mg
Concentration:	1mg/mL
Degree of Substitution: (DOS)	Dye/protein molar ratio is labeled on each vial.
Fluorescence:	Green fluorescence. Excitation/Emission wavelength= 499 nm/523 nm
Storage buffer:	1X phosphate-buffered saline, pH 7.2, with bovine serum albumin and 2 mM sodium azide as preservatives.
Storage:	The dye-protein conjugate is stable for 2~3 months at 4°C. For long-term storage, you can either divide the solution into aliquots and store at -20°C. Avoid multiple thaw-freeze cycles. Or you can add equal volume of glycerol (ACS grade or above) and store the solution at -20°C without aliquoting. The product will be stable for 1 year at -20°C.
Instructions:	The recommended concentration for most immunofluorescent staining is 0.1-10 µg/mL. If the non-specific binding background is high, you may decrease the concentration to minimize the background. You may also centrifuge the conjugate briefly and use the supernatant only for staining. This IgG is highly cross-adsorbed with minimal cross-reaction to human, bovine, horse, rabbit, and swine serum proteins.

For *in vitro* research use only.